

AbFlex® Histone H2A.XS139ph antibody (rAb)

Catalog Nos: 91395, 91396**RRID:** AB_3216368**Application(s):** WB**Reactivity:** Wide Range Predicted**Quantities:** 100 µg, 10 µg**Purification:** Protein A Chromatography**Host:** Mouse**Isotype:** IgG2a**Molecular Weight:** 15 kDa

Background: Histone H2AX phospho Ser139 (H2AX, H2A histone family member X) replaces conventional H2A in a subset of nucleosomes. Nucleosomes wrap and compact DNA into chromatin, limiting DNA accessibility to the cellular machineries that require DNA as a template. Histones thereby play a central role in transcriptional regulation, DNA repair, DNA replication and chromosomal stability. DNA accessibility is regulated via a complex set of post-translational modifications of histones, also called the histone code, and nucleosome remodeling. Histone H2AX is required for checkpoint-mediated arrest of cell cycle progression in response to low doses of ionizing radiation, and for efficient repair of DNA double-strand breaks (DSBs), specifically when modified by C-terminal phosphorylation

Immunogen: This antibody was raised against a phospho-peptide corresponding to phospho-Histone H2AX (Ser139).

Buffer: 140 mM Hepes, pH 7.5, 70 mM NaCl, 32 mM NaOAc, 0.035% sodium azide, and 30% glycerol.

Application Notes:

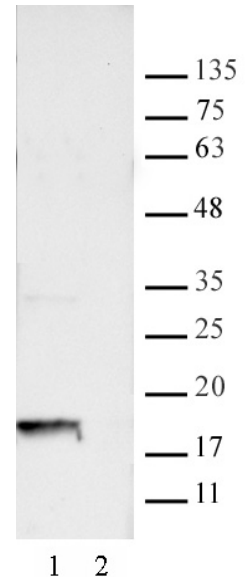
Validated Applications:

WB: 0.5 – 2 µg/ml

*Note: cell lysates prepared by acid extraction is recommended for western blot. Please refer to the method outlines in Active Motif Histone Extraction Kit Manual, Cat. No. 40028, section B, for the protocol we use to prepare acid extracts from cells.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



AbFlex® AbFlex® Histone H2A.XS139ph antibody (rAb) tested by Western blot.

15 µg of HeLa cell nuclear extract was run on SDS-PAGE and probed with AbFlex® Histone H2A.XS139ph antibody at 0.5 µg/ml.